

**PE Mouse Anti-human Drebrin Antibody**  
**\*DBN-N-03, monoclonal\***Catalog number: V103605  
Unit size: 100 tests**Product Details**

---

|                    |                                                                             |
|--------------------|-----------------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving                                                    |
| Concentration      | Lot specific (please consult certificate of analysis for given lot)         |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

---

|                    |            |
|--------------------|------------|
| Species Reactivity | Human      |
| Class              | Primary    |
| Clonality          | Monoclonal |
| Host               | Mouse      |
| Immunogen          | Drebrin    |
| Clone              | DBN-N-03   |
| Conjugate          | PE         |

**Biological Properties**

---

|             |                                                                                                   |
|-------------|---------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions |
| Application | FC (QC TESTED)                                                                                    |

**Applications**

---

This is an antibody that targets Drebrin.